

Bio-Scale DEAE20 Anion Exchange Column

(15 mm x 118 mm) Catalog Number 751-0037

Serial Number: _____
Batch Number: _____
Resin Batch #: _____
Date Tested: _____

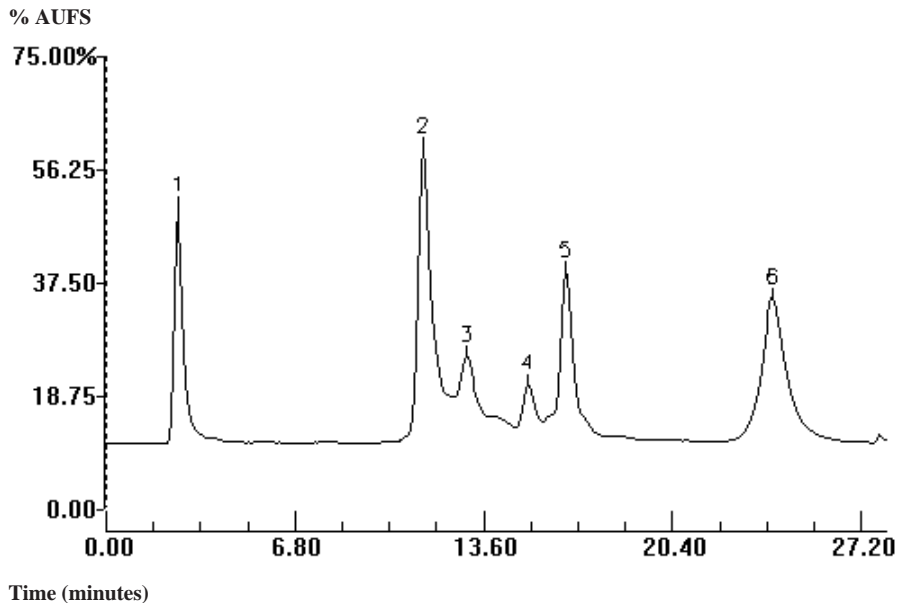
Checked By: _____

Column Specifications

Efficiency (N):	$\geq 12,000$ / meter
Asymmetry (b/a):	0.85 - 1.50
Column Pressure (4.59 ml/min):	< 160 psi

Typical Chromatogram*

Buffer A: 20 mM Tris, pH 7.6	Sample: Protein Standard for Anion Exchange Chromatography (Catalog Number 125-0561)
Buffer B: 20 mM Tris + 1.0 M Sodium Chloride pH 7.6	
Gradient: 0% B in 6 minutes 0 - 50%B in 20 minutes	Inj. Vol: 1000 μ l (9000 μ g Total Protein)
Flow Rate: 4.59 ml/min	Peaks: 1. Myoglobin 2. Conalbumin - A 3. Conalbumin - B 4. Ovalbumin 5. Soybean Trypsin Inhibitor
Detection: UV @ 280 nm, 0.16 AUFS	



* The above chromatogram and test conditions are provided to help resolve any technical problems.